

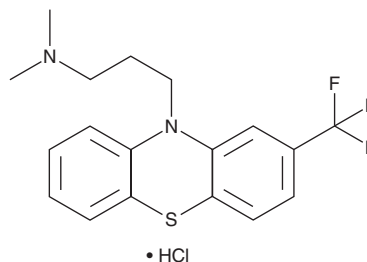
PRODUCT INFORMATION



Triflupromazine (hydrochloride)

Item No. 27610

CAS Registry No.: 1098-60-8
Formal Name: N,N-dimethyl-2-(trifluoromethyl)-10H-phenothiazine-10-propanamine, monohydrochloride
Synonym: Siquil
MF: C₁₈H₁₉F₃N₂S • HCl
FW: 388.9
Purity: ≥98%
UV/Vis.: λ_{max}: 259, 308 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Triflupromazine (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the triflupromazine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Triflupromazine (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of triflupromazine (hydrochloride) in ethanol and DMF is approximately 5 mg/ml and approximately 3 mg/ml in DMSO.

Triflupromazine (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, triflupromazine (hydrochloride) should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Triflupromazine (hydrochloride) has a solubility of approximately 0.09 mg/ml in a 1:10 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Triflupromazine is a phenothiazine with diverse biological activities.¹⁻⁴ It binds to muscarinic receptors in isolated rat corpus striatum (IC₅₀ = 100 μM in a radioligand binding assay).¹ Triflupromazine inhibits serotonin (5-HT) uptake by isolated rat brainstem synaptosomes (IC₅₀ = 0.8 μM).² It inhibits *T. cruzi* infection in mouse peritoneal macrophages when used at a concentration of 12.5 μM.³ Triflupromazine is active against *S. aureus*, shigellae, and vibrios (MICs = 2-100 μg/ml) *in vitro* and is protective against *S. typhimurium* infection in mice when administered at a dose of 30 μg per animal.⁴ Formulations containing triflupromazine were previously used as antipsychotics.

References

1. Snyder, S.H., Greenberg, D., and Yamumura, H.I. Antischizophrenic drugs: Affinity for muscarinic cholinergic receptor sites in the brain predicts extrapyramidal effects. *J. Psychiatr. Res.* **11**, 91-95 (1974).
2. Tuomisto, J. A new modification for studying 5-HT uptake by blood platelets: A re-evaluation of tricyclic antidepressants as uptake inhibitors. *J. Pharm. Pharmacol.* **26**(2), 92-100 (1974).
3. De Castro, S.L., Soeiro, M.N., and De Meirelles Mde, N. *Trypanosoma cruzi*: Effect of phenothiazines on the parasite and its interaction with host cells. *Mem. Inst. Oswaldo Cruz* **87**(2), 209-215 (1992).
4. Dastidar, S.G., Debnath, S., Mazumdar, K., et al. Triflupromazine: A microbicide non-antibiotic compound. *Acta Microbiol. Immuno. Hung.* **51**(1-2), 75-83 (2004).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/09/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM